

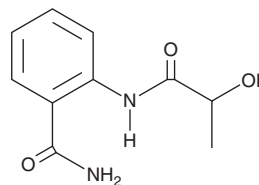
PRODUCT INFORMATION



2-((2-Hydroxypropionyl)amino)benzamide

Item No. 40915

CAS Registry No.: 18326-27-7
Formal Name: 2-[(2-hydroxy-1-oxopropyl)amino]-benzamide
Synonyms: N-(2-hydroxypropanoyl)-2-Aminobenzoic Acid Amide
MF: C₁₀H₁₂N₂O₃
FW: 208.2
Purity: ≥70%
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years
Item Origin: Bacterium/*Penicillium atrovenetum*



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

2-((2-Hydroxypropionyl)amino)benzamide is supplied as a solid. A stock solution may be made by dissolving the 2-((2-hydroxypropionyl)amino)benzamide in the solvent of choice, which should be purged with an inert gas. 2-((2-Hydroxypropionyl)amino)benzamide is soluble in ethanol, methanol, dichloromethane, and DMSO.

Description

2-((2-Hydroxypropionyl)amino)benzamide is a fungal metabolite that has been found in *Penicillium* and has angiogenic activity.¹ It induces angiogenesis in a zebrafish larval model of vascular insufficiency induced by the VEGFR inhibitor PTK787 (vatalanib; Item No. 14868) when added to the tank water at a concentration of 100 µg/ml.

Reference

1. Xie, D., Chen, H., et al. Secondary metabolites from marine derived fungus *Penicillium chrysogenum* Y19-1 with proangiogenic and antithrombotic activities. *Biochem. Syst. Ecol.* **107**, 104625 (2023).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 05/16/2024

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM